

ANGPTL3 polyclonal antibody

This antibody is covered by our [Worry-Free Guarantee](#).

Ordering Information

[Order Online »](#)

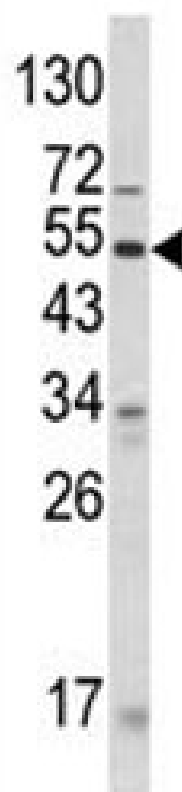
ENZ-ABS868-0400	0.4ml
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of ANGPTL3 antibody and mouse liver tissue lysate. Predicted molecular weight 50 ~ 63 kDa depending on glycosylation level.



Western blot analysis of ANGPTL3 antibody and human CEM lysate. Predicted molecular weight 50 ~ 63 kDa depending on glycosylation level.

Handling & Storage

Handling	Aliquot the ANGPTL3 antibody and store frozen at -20°C or colder. Avoid freeze/thaw cycles.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	Angiopoietin-related protein 3, ANG-5, Angiopoietin-5
Application	ELISA, WB
Formulation	Liquid. In 1X PBS, pH 7.4, containing 0.09% sodium azide.
Host	Rabbit
Immunogen	Human ANGPTL3 (aa 82-114).
Isotype	Ig
Purity Detail	Purified.
Recommendation Dilutions/Conditions	Western blot (1:1,000)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Human, Mouse
UniProt ID	Q9Y5C1
Worry-free Guarantee	This antibody is covered by our Worry-Free Guarantee .



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)